

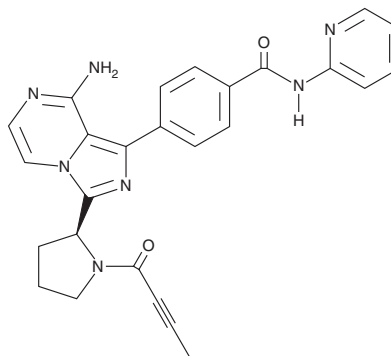
PRODUCT INFORMATION



Acalabrutinib

Item No. 19899

CAS Registry No.: 1420477-60-6
Formal Name: 4-[8-amino-3-[(2S)-1-(1-oxo-2-butyn-1-yl)-2-pyrrolidinyl]imidazo[1,5-a]pyrazin-1-yl]-N-2-pyridinyl-benzamide
Synonym: ACP-196
MF: C₂₆H₂₃N₇O₂
FW: 465.5
Purity: ≥95%
UV/Vis.: λ_{max}: 229, 283, 336 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Acalabrutinib is supplied as a crystalline solid. A stock solution may be made by dissolving the acalabrutinib in the solvent of choice, which should be purged with an inert gas. Acalabrutinib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of acalabrutinib in ethanol is approximately 15 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Acalabrutinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, acalabrutinib should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Acalabrutinib has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Acalabrutinib is an irreversible inhibitor of Bruton's tyrosine kinase (BTK). It can dose-dependently inhibit B cell receptor signaling in primary chronic lymphocytic leukemia cells.¹ In kinase-inhibition assays, acalabrutinib was more selective than ibrutinib (Item No. 16274) for inhibiting BTK and did not inhibit EGFR, Itk, or Txk signaling.¹

Reference

1. Burd, J.C., Harrington, B., O'Brien, S., *et al.* Acalabrutinib (ACP-196) in relapsed chronic lymphocytic leukemia. *N. Engl. J. Med.* **374**(4), 323-332 (2016).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/09/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM